

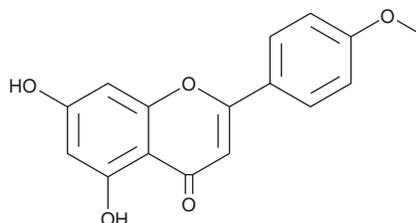
PRODUCT INFORMATION



Acacetin

Item No. 20827

CAS Registry No.: 480-44-4
Formal Name: 5,7-dihydroxy-2-(4-methoxyphenyl)-4H-1-benzopyran-4-one
Synonyms: LY064233, NSC 76061
MF: C₁₆H₁₂O₅
FW: 284.3
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 268, 327 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Acacetin is supplied as a crystalline solid. A stock solution may be made by dissolving the acacetin in the solvent of choice. Acacetin is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of acacetin in these solvents is approximately 15 mg/ml.

Acacetin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, acacetin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Acacetin has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Acacetin is an O-methylated flavone found in various plants. It is reported to demonstrate spasmolytic, antinociceptive, anti-inflammatory, and antioxidant activity in various research models.^{1,2}

References

- Zielińska, S. and Matkowski, A. Phytochemistry and bioactivity of aromatic and medicinal plants from the genus *Agastache* (*Lamiaceae*). *Phytochem. Rev.* **13**(2), 391-416 (2014).
- Jaganathan, S.K. and Mandal, M. Antiproliferative effects of honey and of its polyphenols: A review. *J. Biomed. Biotechnol.* **2009**, 830616 (2009).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/26/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM